



**CERTIFICATE OF ANALYSIS**

**Product Description:** A $\beta$ 1-42 Peptide (4G8)  
**Catalog Number:** C003A-2  
**Lot Number:** A0080393  
**Expiration Date:** 30 SEP 2026  
**Storage Temperature:**  $\leq -70^{\circ}\text{C}$

**Concentration Assignment:**

Analyte	Stock Concentration	Unit of Measure
A $\beta$ 42	80400	pg/mL

**Dilution to Top of Curve:** 40 X

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		13 MAY 2024